

Macrophage marker monoclonal antibody (HAM56) (Ready-to- Use)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-30935	~6 µg - 6ml
-----------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	IHC (PS)
Clone	HAM56
Crossreactivity	Cross-reacts with a subpopulation of endothelial cells including those of capillaries and smaller blood vessels.
Formulation	Liquid. <i>Ready-to-use</i> solution in PBS containing 0.1% sodium azide and 1% BSA.
Host	Mouse
Immunogen	Sonicated human alveolar macrophages.
Isotype	IgM
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections; this antibody can be used for staining of paraffin-embedded tissue sections fixed in neutral buffered formalin and is compatible with most commonly used histological fixatives. Choose a good quality enzyme-linked anti-mouse IgM antibody for detection). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human, Monkey
Specificity	Recognizes fixed macrophages such as “tingible macrophages”, interdigitating macrophages of lymph nodes or tissue macrophages.

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)